

# OVA 323-339

Chicken egg albumin peptide

Catalog code: vac-isq

<https://www.invivogen.com/ova-peptide>

For research use only

Version 22F27-MM

## PRODUCT INFORMATION

### Contents

- 1 mg OVA 323-339 provided lyophilized

### Storage and stability

- OVA 323-339 is shipped at room temperature. Upon receipt, store at -20°C.

*Note: OVA 323-339 is hygroscopic and light sensitive. Keep tightly closed and protect from light.*

- Upon resuspension, prepare aliquots of OVA 323-339 and store at -20°C. Resuspended product is stable for 6 months when properly stored. Avoid repeated freeze-thaw cycles.

### Quality Control

- Purity: ≥ 95% (UHPLC)
- The absence of bacterial contamination (e.g. lipoproteins and endotoxins) has been confirmed using HEK-Blue™ TLR2 and HEK-Blue™ TLR4 cells.

## DESCRIPTION

OVA 323-339 encompasses an allergenic and antigenic epitope of the ovalbumin protein<sup>1</sup>. OVA 323-339 peptide binds to I-A(d) Major Histocompatibility Complex (MHC) class II protein<sup>2</sup>. OVA 323-339 has been used extensively to study the nature of class II MHC-peptide binding and T-cell activation<sup>3-4</sup>.

**Sequence:** ISQAVHAAHAEINEAGR

(H - Ile - Ser - Gln - Ala - Val - His - Ala - Ala - His - Ala - Glu - Ile - Asn - Glu - Ala - Gly - Arg - OH)

1. **McFarland B.J. et al. 1999.** Ovalbumin(323-339) peptide binds to the major histocompatibility complex class II I-A(d) protein using two functionally distinct registers. *Biochemistry*, 38(50):16663-70. 2. **Johnsen G & Elsayed S. 1990.** Antigenic and allergenic determinants of ovalbumin-111. MHC Ia-binding peptide (OA 323-339) interacts with human and rabbit specific antibodies. *Mol Immunol*, 27:821-7. 3. **Bickert T. et al., 2007.** Murine polyomavirus-like particles induce maturation of bone marrow-derived dendritic cells and proliferation of T cells. *Med Microbiol Immunol* 196:31-9. 4. **Szymczak-Workman A.L. et al., 2009.** Cutting edge: regulatory T cells do not require stimulation through their TCR to suppress. *J Immunol* 1:182.

## CHEMICAL PROPERTIES

Synonym: ISQAVHAAHAEINEAGR

Chemical formula: C<sub>74</sub>H<sub>120</sub>N<sub>26</sub>O<sub>25</sub>

Molecular weight: 1773.9 g/mol

Solubility: 1 mg/ml in water

## METHODS

Preparation of stock solution (1 mg/ml)

1. Add 1 ml water to 1 mg of OVA 323-339.
2. Mix the solution by pipetting up and down.

## APPLICATIONS

OVA 323-339 can be used for Enzyme-Linked ImmunoSpot (ELISPOT) assay at a working concentration of 10 µg/ml.

Other applications have not been tested.

## RELATED PRODUCTS

Product	Description	Cat.Code
EndoFit™ Ovalbumin	Ova for <i>in vivo</i> use	vac-pova
Ovabumin	Ova for detection use	vac-stova
OVA 257-264	For detection; ELISPOT	vac-sin

## TECHNICAL SUPPORT

InvivoGen USA (Toll-Free): 888-457-5873

InvivoGen USA (International): +1 (858) 457-5873

InvivoGen Europe: +33 (0) 5-62-71-69-39

InvivoGen Asia: +852 3622-3480

E-mail: [info@invivogen.com](mailto:info@invivogen.com)